

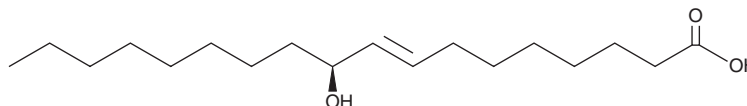
# PRODUCT INFORMATION



## 10(S)-HOME

Item No. 33246

**CAS Registry No.:** 263023-34-3  
**Formal Name:** 10S-hydroxy-8E-octadecenoic acid  
**Synonym:** HOD  
**MF:** C<sub>18</sub>H<sub>34</sub>O<sub>3</sub>  
**FW:** 298.5  
**Purity:** ≥98%  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

10(S)-HOME is supplied as a crystalline solid. A stock solution may be made by dissolving the 10(S)-HOME in the solvent of choice, which should be purged with an inert gas. 10(S)-HOME is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 10(S)-HOME in these solvents is approximately 100, 20, and 15 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of 10(S)-HOME can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 10(S)-HOME in PBS (pH 7.2) is approximately 0.25 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

10(S)-HOME is an oxylipin. It is an intermediate in the conversion of oleic acid (Item Nos. 90260 | 24659) to the antibacterial hydroxy fatty acid 7,10-dihydroxy-8(E)-octadecenoic acid (Item No. 33447) in *P. aeruginosa*.<sup>1</sup>

### Reference

1. Kim, H., Gardner, H.W., and Hou, C.T. 10(S)-Hydroxy-8(E)-octadecenoic acid, an intermediate in the conversion of oleic acid to 7,10-dihydroxy-8(E)-octadecenoic acid. *J. Am. Oil Chem. Soc.* **77(1)**, 95-99 (2000).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 03/01/2021

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897

[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM